

## POLD3 Antibody [#C15428]

<b>Catalog Number:</b>	C15428
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q15054
<b>Other Names:</b>	DNA polymerase subunit delta-3; DNA polymerase subunit delta p66
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without $Mg^{2+}$ and $Ca^{2+}$ ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from internal of human POLD3.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	POLD3 antibody detects endogenous levels of total POLD3 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:40000
<b>References:</b>	Nomura N., DNA Res. 1:27-35(1994). Ota T., Nat. Genet. 36:40-45(2004). Mural R.J., Submitted (JUL-2005) to the EMBL/GenBank/DDBJ databases.